

BRD4 NanoBRET TE Intracellular Assay

BRD4 [NanoBRET Target Engagement Intracellular Assay](#) for screening of compounds binding to the target, competitive binding assay format

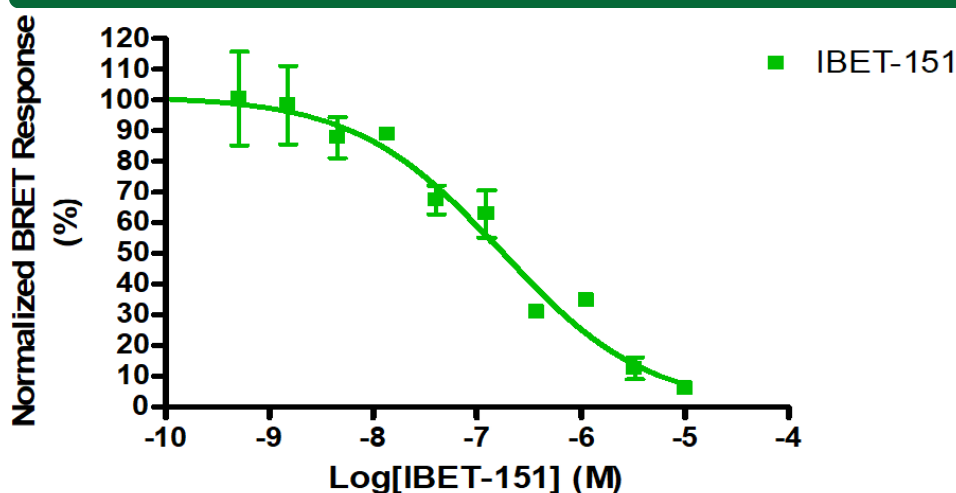
Target Overview

Target	BRD4
HGNC Symbol	BRD4
Synonyms	HUNK1; MCAP; CAP; HUNK1
Cell Line	HEK293 human epithelial cell line from ATCC

Assay Properties

Assay Format	384-well plate format using Promega's NanoBRET Intracellular Target Engagement Assay
Assay Protocol	HEK293 cells transiently expressing NanoLuc-BRD4 Fusion Vector were seeded into the wells of 384-well plates. The cells were pre-treated with the NanoBRET BRD Tracer and then treated with reference compound IBET-151 for 1 hour. The BRET signal was measured on an Envision 2104 Multilabel Reader. IC50 value was calculated and IC50 curve was plotted using the GraphPad Prism program based on a sigmoidal dose response equation.
Readout	Binding of compound to target
Reference Compound IC50s	Compounds IC50 (M) IBET-151 1.710e-007
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website Cell-based Epigenetic Assays

Reference compound IC50 for BRD4



	IBET-151 (M)
HILLSLOPE	-0.6237
EC50	1.710e-007

Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/brd4_nano_malvern

